



Product Data Sheet

Product Name:	Charybdotoxin	
Catalog Number:	ØJÈ28244 (0.1 mg)	Lot Number: See label on vial
Sequence:	Pyr-Phe-Thr-Asn-Val-Ser-Cys-Thr-Thr-Ser-Lys-Glu-Cys-Trp-Ser-Val-Cys-Gln-Arg-Leu-His-Asn-Thr-Ser-Arg-Gly-Lys-Cys-Met-Asn-Lys-Lys-Cys-Arg-Cys-Tyr-Ser-OH (Disulfide bridge: 7-28, 13-33 and 17-35) (3-letter code) Pyr-FTNVSCTTSKECWSVCQRLHNTSRGKCMNKKCRCYS (Disulfide bridge: 7-28, 13-33 and 17-35) (1-letter code)	
Molecular Weight:	4296.0	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Charybdotoxin peptide is freely soluble in water.

Storage: Charybdotoxin peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description: Charybdotoxin (ChTX) is a Ca^{2+} -activated K^{+} channel blocker. It depolarizes peripheral T lymphocytes and blocks their mitogen-induced proliferation. ChTX is a highly basic peptide isolated from venom of the scorpion, *Leiurus quinquestriatus hebraeus*. Ref: Leonard, R. et al. *Proc. Natl. Acad. Sci. USA* **89**, 10094 (1992); Gimenez-Gallego, *Proc. Natl. Acad. Sci. USA* **85**, 3329 (1988). Sugg, E. et al. *J. Biol. Chem.* 265, 18745 (1990).

Published Citations:

- Nemargut, JP. et al. *Vision Research* **49**, 388 (2009).
- Bayram, Z. et al. *Regulatory Peptides* doi:10.1016/j.regpep.2009.12.023 (2009).

Related Products:

Name	Cat #	Size
Iberitoxin (IbTX)	AS-60763	0.1 mg
Pyr-FTDVDCSVSKECWSVCKDLFGVDRGKCMGKKRCYQ (Disulfide bridge: C7-C28,C13-C33,C17-C35)		

For Research Use Only